## **MB Acid**

MB (methylene blue) acid can be conjugated to peptides, proteins, drugs, polymeric materials and biomolecules with primary amine groups. The conjugate will have a blue color and be able to complex with nucleic acids.



## **Product attributes**

## **Product Description**

Methylene Blue (MB) is a commonly used redox indicator in nucleic acid research. It is also studied for its use in medical applications as well as being used as a general biological stain. Reactive formats of MB can be conjugated to peptides, proteins, drugs, polymeric materials and other biomolecules. The conjugate will have a blue color and be able to complex with nucleic acids.

- MW: 392
- Store at 4°C and protected from light

We also offer a selection of chemically reactive formats for use in labeling biomolecules such as proteins and nucleic acids. See the table below for our full list of metheylene blue derivatives and formats.

## **Methylene Blue Derivatives**

Product	Size	Catalog No.	Features
MB Acid	5 mg	<u>40076</u>	Free acid form
MB Succinimidyl Ester	5 mg	40075	Amine-reactive chemistry for labeling proteins
MB-Maleimide	1 mg	40118	Thiol-reactive chemistry for labeling proteins
MB-DBCO	1 mg	<u>40114</u>	Allows bioorthogonal conjugation to label azide containing molecules
MB-Methyltetrazine	1 mg	<u>40115</u>	Allows labeling of TCO tagged molecules
MB-TCO	1 mg	40116	Allows labeling of tetrazine tagged molecules
MB-Azide	1 mg	40117	Allows labeling alkyne, BCN, or phosphine-containing molecules.

See our other reactive DNA/RNA binding dyes.

This datasheet was generated on November 10, 2025 at 08:07:37 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/mb-acid/">https://biotium.com/product/mb-acid/</a>